



Filling MicroClime Environmental Lids

Introduction

The MicroClime Environmental lid minimizes edge effects and preserves the concentrations of solutions in microplate wells. Unlike standard microplate lids, the MicroClime lid incorporates a unique, fluid-absorbing matrix. The absorbed fluid creates a vapor barrier that protects the sample from the exterior environment, especially near the edges of the microplate.

The MicroClime lid is easy to use—simply place it on or lift it off a microplate. No specialized tools are required, and it is compatible with robotic plate lidding systems. The lid can be filled with many different fluids, using either a hand-held syringe or automated bulk fillers.

When filled with anhydrous DMSO, the MicroClime lid prevents evaporation and hydration of DMSO in compound library plates during storage. It also stops evaporation of ultra-low-volume DMSO droplets in screening assays. The lid can also be filled with water to prevent evaporation in aqueous assays.

